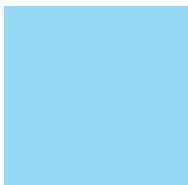


**Mini-PROTEAN® Tetra Cell**



**The Best Just Got Better**





The Best Just Got Better  
**Mini-PROTEAN Tetra Cell**

For over 20 years, the Mini-PROTEAN® system has set the standard for excellence in protein electrophoresis. The new Mini-PROTEAN Tetra cell is the next-generation mini cell for 1-D vertical gel electrophoresis. Easier to assemble, leakproof, and with a four-gel capacity that doubles the output of the Mini-PROTEAN family, the Mini-PROTEAN Tetra cell combines traditional Mini-PROTEAN features with higher throughput and even greater reliability.

For more information on Bio-Rad's vertical electrophoresis cells, refer to the 2006/07 Life Science Research Product catalog, or go to [www.bio-rad.com/verticalelectro/](http://www.bio-rad.com/verticalelectro/)

The Mini-PROTEAN Tetra cell is ideal for your vertical mini gel electrophoresis needs. Easy to use, easy to assemble, robust, and versatile, the Tetra cell accommodates up to four handcast or Ready Gel® precast gels. Casting and running gels have never been quicker or easier.

- Run up to 4 mini gels in less than an hour
- Run four 2-D gels in less than a day
- Evaluate sample preparation conditions
- Screen new samples and process discovery projects



### Casting Advantages

- Casting frames with simple cam closure provide precision alignment on any flat surface
- Patented side-by-side casting stand\* allows access to two gels simultaneously, and the spring-loaded lever creates a tight seal against the thermoplastic gaskets
- Thermoplastic casting gaskets form a tight seal with the glass plates to ensure leak-free casting
- Ground-glass plates with permanently bonded spacers guarantee perfect alignment and leak-free casting
- Thick glass spacer plates reduce breakage
- Glass plates and combs are labeled with thickness and number of wells for instant identification
- Patented plastic combs\* do not inhibit polymerization and have a built-in ridge to eliminate air contact during gel casting for uniform gel polymerization

### Loading and Running Innovations

- Patented sample loading guides\*\* prevent skipped or repeated loading lanes
- Cell runs 1–4 gels using 2 running modules

\* US patent 6,162,342.

\*\* US patent 5,656,145.

### Modular Cells for Many Applications

Interchangeable modules easily convert a Mini-PROTEAN Tetra cell from one application to another. Each module fits into the same buffer tank and lid to form a complete cell.

The modular configurations are:

- Mini-PROTEAN Tetra electrophoresis cell — for running handcast or precast polyacrylamide gel electrophoresis (PAGE) and sodium dodecyl sulfate (SDS)-PAGE gels
- Mini Trans-Blot® electrophoretic transfer cell — for western blotting

### Configure Your Own Electrophoresis Cell

To configure your own electrophoresis cell, order the Mini-PROTEAN Tetra cell (catalog #165-8004) and select a casting module from the ordering information on the next page. For a preset configuration, order catalog #165-8000 (10-well, 0.75 mm) or #165-8001 (10-well, 1.0 mm).

### Specifications

Number of gels	1–4
Precast gels	Ready Gel precast gels
Handcast gels	Cast using Mini-PROTEAN spacer plates
Glass plate size (W x L)	
Short plate	10.1 x 7.3 cm
Spacer plate	10.1 x 8.2 cm
Gel size (W x L)	
Handcast	8.3 x 7.3 cm
Precast	8.6 x 6.8 cm
Total buffer volume for 2 gels	700 ml
Total buffer volume for 4 gels	1,000 ml
Typical run times for SDS-PAGE	35–45 min (at 200 V constant)
Recommended power supply	PowerPac™ Basic or PowerPac™ HC
Dimensions (W x L x H)	12 x 16 x 18 cm
Weight	1.0 kg



## Ordering Information

Catalog #	Description	Description	0.75 mm	1.0 mm	1.5 mm	
<b>Mini-PROTEAN Tetra Cells</b>						
165-8000	<b>Mini-PROTEAN Tetra Cell</b> , 10-well, 0.75 mm thickness; 4-gel system includes 5 combs, 5 sets of glass plates, 2 casting stands, 4 casting frames, sample loading guide, electrode assembly, companion running module, tank, lid with power cables, mini cell buffer dam	<b>Casting Modules**</b>				
165-8001	<b>Mini-PROTEAN Tetra Cell</b> , 10-well, 1.0 mm thickness; 4-gel system includes 5 combs, 5 sets of glass plates, 2 casting stands, 4 casting frames, sample loading guide, electrode assembly, companion running module, tank, lid with power cables, mini cell buffer dam	5-Well	165-8008	165-8013	165-8019	
165-8002*	<b>Mini-PROTEAN Tetra Cell</b> , 10-well, 0.75 mm thickness; 2-gel system includes 5 combs, 5 sets of glass plates, casting stand, 2 casting frames, sample loading guide, electrode assembly, tank, lid with power cables, mini cell buffer dam	9-Well	165-8009	165-8014	165-8020	
165-8003*	<b>Mini-PROTEAN Tetra Cell</b> , 10-well, 1.0 mm thickness; 2-gel system includes 5 combs, 5 sets of glass plates, casting stand, 2 casting frames, sample loading guide, electrode assembly, tank, lid with power cables, mini cell buffer dam	10-Well	165-8010	165-8015	165-8021	
165-8004	<b>Mini-PROTEAN Tetra Cell for Ready Gel Precast Gels</b> , 4-gel system includes electrode assembly, companion running module, tank, lid with power cables, mini cell buffer dam	15-Well	165-8011	165-8016	165-8022	
		Prep/2-D Well	165-8012	165-8017	165-8023	
		IPG Well	—	165-8018	165-8024	
		Catalog #	Description			
<b>Mini-PROTEAN Tetra Systems</b>						
165-8025	<b>Mini-PROTEAN Tetra Cell and PowerPac Basic Power Supply</b> , includes 165-8001 and 164-5050	<b>Handcast Gel Accessories and Replacement Parts</b>				
165-8026	<b>Mini-PROTEAN Tetra Cell and PowerPac™ Universal Power Supply</b> , includes 165-8001 and 164-5070	165-3303	<b>Mini-PROTEAN Casting Stand</b>			
165-8027	<b>Mini-PROTEAN Tetra Cell and PowerPac HC Power Supply</b> , includes 165-8001 and 164-5052	165-3304	<b>Mini-PROTEAN Casting Frame</b>			
165-8028	<b>Mini-PROTEAN Tetra Cell and PowerPac™ HV Power Supply</b> , includes 165-8001 and 164-5056	165-3305	<b>Mini-PROTEAN Casting Stand Gaskets</b> , replacement, 2			
165-8029	<b>Mini-PROTEAN Tetra Cell and Mini Trans-Blot Module</b> , includes 165-8001 and 170-3935	165-3308	<b>Short Plates</b> , 5			
165-8030	<b>Mini-PROTEAN Tetra Cell for Ready Gel Precast Gels and Mini Trans-Blot Module</b> , includes 165-8004 and 170-3935	165-3309	<b>Spacer Plates With 0.5 mm Integrated Spacers</b> , 5			
165-8033	<b>Mini-PROTEAN Tetra Cell, Mini Trans-Blot Module, and PowerPac Basic Power Supply</b> , includes 165-8001, 170-3935, and 164-5050	165-3310	<b>Spacer Plates With 0.75 mm Integrated Spacers</b> , 5			
165-8034	<b>Mini-PROTEAN Tetra Cell for Ready Gel Precast Gels, Mini Trans-Blot Module, and PowerPac Basic Power Supply</b> , includes 165-8004, 170-3935, and 164-5050	165-3311	<b>Spacer Plates With 1.0 mm Integrated Spacers</b> , 5			
165-8035	<b>Mini-PROTEAN Tetra Cell, Mini Trans-Blot Module, and PowerPac HC Power Supply</b> , includes 165-8001, 170-3935, and 164-5052	165-3312	<b>Spacer Plates With 1.5 mm Integrated Spacers</b> , 5			
165-8036	<b>Mini-PROTEAN Tetra Cell for Ready Gel Precast Gels, Mini Trans-Blot Module, and PowerPac HC Power Supply</b> , includes 165-8004, 170-3935, and 164-5052		Description	0.75 mm	1.0 mm	1.5 mm
			<b>Combs</b>			
			5-Well	165-3352	165-3357	165-3363
			9-Well	165-3353	165-3358	165-3364
			10-Well	165-3354	165-3359	165-3365
			15-Well	165-3355	165-3360	165-3366
			Prep/2-D Well	165-3356	165-3361	165-3367
			IPG Well	—	165-3362	165-3368
			Catalog #	Description		
			<b>Other Replacement Parts</b>			
			165-8037	<b>Mini-PROTEAN Tetra Electrode Assembly</b>		
			165-8038	<b>Mini-PROTEAN Tetra Companion Running Module Buffer Tank</b> , replacement		
			165-8039	<b>Buffer Tank and Lid</b> , replacement		
			165-8040	<b>Cell Lid With Power Cables</b>		
			165-8041	<b>Sample Loading Guide</b> , 9-well (red)		
			165-3201	<b>Sample Loading Guide</b> , 10-well (yellow)		
			165-3146	<b>Sample Loading Guide</b> , 12-well (green)		
			165-3203	<b>Sample Loading Guide</b> , 15-well (blue)		
			165-3132	<b>Mini Cell Buffer Dams</b> , 2		
			165-3130	<b>Replacement Gaskets</b> , for electrophoresis assembly, green, 2		
			165-3149	<b>Empty Ready Gel Cassettes</b> , 10		
			161-0990			

\* The 2-gel systems do not include the companion running module.

\*\* Each casting module includes 2 casting stands, 4 casting frames, 5 combs, 5 sets of plates, and the appropriate sample loading guide.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400 Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65 Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300 Israel 03 963 6050 Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500 Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723 Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000